

OCFD03-FT4-K

Protocole/Protocol : 45mn

Lot N°/Batch # : 376B

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-OCFD03-FT4	376A	27/07/2022
CT-OCFD03-FT4	147B	25/01/2023
CAL0 OCFD03-FT4-K	20F	03/10/2022
CAL1 OCFD03-FT4-K	20F	03/10/2022
CAL2 OCFD03-FT4-K	20F	03/10/2022
CAL3 OCFD03-FT4-K	20F	03/10/2022
CAL4 OCFD03-FT4-K	20F	03/10/2022
CAL5 OCFD03-FT4-K	20F	03/10/2022
CAL6 OCFD03-FT4-K	20F	03/10/2022
CONT OCFD03-FT4-K	20F	03/10/2022
OCFD03-FT4-K MO	002	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	pg/mL	B0/T %	50.73
CAL.1	3.00	pg/mL	B/B0	77.74
CAL.2	6.00	pg/mL	B/B0	65.03
CAL.3	10.50	pg/mL	B/B0	53.18
CAL.4	20.00	pg/mL	B/B0	37.71
CAL.5	42.00	pg/mL	B/B0	19.87
CAL.6	68.00	pg/mL	B/B0	11.11

Activité totale / Total activity : 75436 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	pg/mL	7.36 - 11.04	8.65

Date du contrôle
Control Date

17/05/2022

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

27/07/2022

